

PRODUCT INFORMATION



BI-907828

Item No. 44353

CAS Registry No.: 2095116-40-6

Formal Name: (2'S,3'S, 3'aS,10'aS)-6-chloro-3'-(3-chloro-2-fluorophenyl)-1'-(cyclopropylmethyl)-1,2,3',3'a,10',10'a-hexahydro-6'-methyl-2-oxo-spiro[3H-indole-3,2'(1'H)-pyrrolo[2',3':4,5]pyrrolo[1,2-b]indazole]-7'carboxylic acid

Synonym: Brigimadlin

MF: C₃₁H₂₅Cl₂FN₄O₃

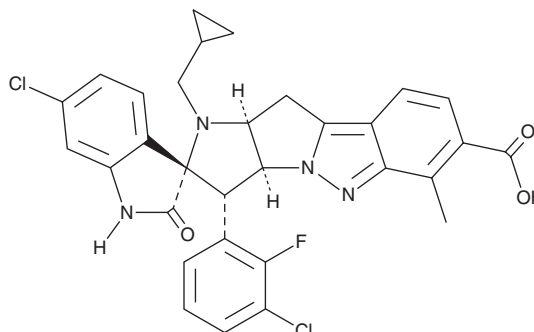
FW: 591.5

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

BI-907828 is supplied as a solid. A stock solution may be made by dissolving the BI-907828 in the solvent of choice, which should be purged with an inert gas. BI-907828 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

BI-907828 is an inhibitor of the protein-protein interaction between p53 and murine double minute 2 (MDM2; IC₅₀ = 2 nM in a cell-free assay).¹ It is selective for the p53-MDM2 interaction over the p53-MDMX interaction (IC₅₀ = 7,944 nM). BI-907828 selectively inhibits the proliferation of Colon-26 colon cancer cells, which express wild-type p53, over 4T1 breast and MC-38 colon cancer cells, which express mutant p53 (IC₅₀s = 259, 5,579, and 9,782 nM, respectively). It inhibits the proliferation of SJSA-1 osteosarcoma cells, BJ human foreskin fibroblasts, and hTERT RPE-1 retinal pigment epithelial cells but only induces apoptosis in SJSA-1 cells. BI-907828 induces tumor regression in an SJSA-1 mouse xenograft model and patient-derived xenograft (PDX) models.

Reference

1. Gollner, A., Rudolph, D., Weyer-Czernilofsky, U., *et al.* Discovery and characterization of brigimadlin, a novel and highly potent MDM2-p53 antagonist suitable for intermittent dose schedules. *Mol. Cancer Ther.* **23**(12), 1689-1702 (2024).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/20/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM